

Validation data for ODN 2006 Biotin

<https://www.invivogen.com/odn2006-biotin>

For research use only

Version 24E06-AK

ODN 2006 Biotin is a Class B CpG oligonucleotide (ODN). It is a ligand of choice for human Toll-like receptor 9 (TLR9). Activation of TLR9 with ODN 2006 triggers NF- κ B- and interferon regulatory factor (IRF)-mediated pro-inflammatory responses. The labeled ligand retains the biological activity of ODN 2006. It can be used to evaluate CpG ODN binding, cellular uptake, and localization by microscopy [data not shown] or flow cytometry (**Figure 1**).

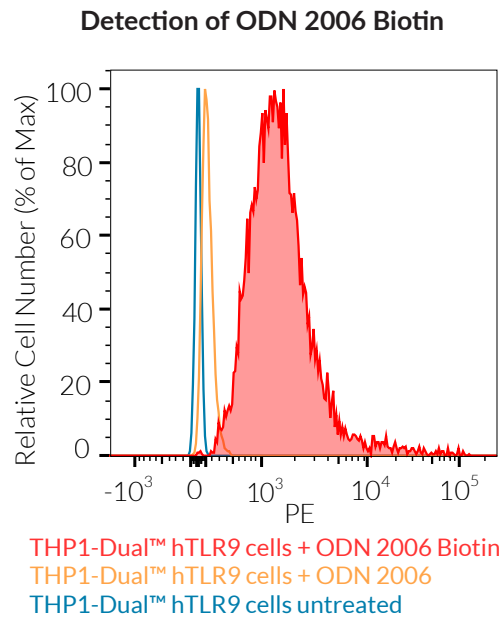


Figure 1. Intracellular detection of ODN 2006 Biotin. THP1-Dual™ hTLR9 cells were incubated with ODN 2006 unmodified or labeled with Biotin for 16h - 24h. After, the cells were fixed and permeabilized, following the incubation with an Anti-Biotin-PE secondary antibody. The intracellular staining was analyzed by flow cytometry.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Asia: +852 3622-3480
E-mail: info@invivogen.com